

# (D-Pro1,2)-retro-(D-HIV-TAT(48-57))

## Jnk inhibitor

Useful as a control peptide in studies using D-JNKi1 (BML-EI355). This peptide is the cell transporter, and does not contain the JIP (JNK inhibitor peptide) sequence.

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Appearance</b>	Solid.
<b>Formulation</b>	Lyophilized.
<b>MW</b>	1589.5
<b>Purity</b>	≥98% (TLC)
<b>Sequence</b>	H-DPro-DPro-DArg-DArg-DArg-DGln-DArg-DArg-DLys-DLys-DArg-Gly-NH <sub>2</sub>
<b>Solubility</b>	Soluble in Water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)